



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-08000

Product Description: Rat BNP Calibrator
Catalog Number: C03KF-2
Lot Number: A0040187
Expiration Date: 30 JUN 2019
Storage Temperature: ≤ -70°C

Stock Concentration of Materials Tested:

	Assay
	Rat BNP
Calibrator (pg/mL)	41200

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 MAR 2017